
LSD1 Protein

(Catalog No. E24027)

Description

Human LSD1 [also known as KDM1 (lysine (K)-specific demethylase 1) or AOF2, amine oxidase (flavin containing) domain 2] GenBank Accession No. NM_015013, amino acids 158-end with N-terminal GST tag.

Molecular Weight (kDa)

103 kDa

Formulation

50 mM Sodium phosphate buffer (pH 7.4)

Source

Expressed in E. coli expression system

Size

50 µg

Storage Temp & Conditions

>6 months at -80°C

Purification Notes

Purity>40%

Application Notes

Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Ordering Information:

Products	Size	Cat. No.
LSD1 Protein	50 µg	E24027 -1

This product is for research purposes only. Not intended for use in diagnostic procedures.